

Product Number:	103187
Lot Number:	502602
Expiration:	29-Apr-2021

Component	Item Number	Lot Number
RQT Plate	103095	030304
RGP Reagent	103159	N/A ¹

¹ RGP is not Kit Lot Specific

Means of Analysis	Result
Simoa SR-X Analyzer	PASS

Review/Approval

Shivani Goel
Name

QC Manager
Title

Shivani Goel 09 NOV 2020
Signature/Date